

Product Information



Prostaglandin Screening Library III (96-Well)

Item No. 10503 • Batch No. 0469449

The composition of the library will always vary somewhat depending upon our inventory.

Page 1 of 3

Plate	Well	Contents	Item Number
1	A1	Unused	
1	A2	Prostaglandin A ₁	10010
1	A3	Prostaglandin A ₂	10210
1	A4	BW 245C	12050
1	A5	1a,1b-dihomo Prostaglandin F _{2α}	16050
1	A6	11-deoxy Prostaglandin F _{2α}	16500
1	A7	6α-Prostaglandin I ₁	18110
1	A8	Prostaglandin J ₂	18500
1	A9	Tafluprost (free acid)	10005439
1	A10	8-iso-15-keto Prostaglandin F _{2β}	10008539
1	A11	Unused	
1	A12	Unused	
1	B1	Unused	
1	B2	Prostaglandin A ₁ ethyl ester	10020
1	B3	13,14-dihydro-15-keto Prostaglandin A ₂	10260
1	B4	BW 246C	12055
1	B5	Prostaglandin F _{2α} 1,11-lactone	16110
1	B6	11-deoxy Prostaglandin F _{2β}	16510
1	B7	6β-Prostaglandin I ₁	18120
1	B8	15-deoxy-Δ ^{12,14} -Prostaglandin J ₂	18570
1	B9	Tafluprost	10005440
1	B10	13,14-dihydro-15-keto Prostaglandin J ₂	10008840
1	B11	Unused	
1	B12	Unused	
1	C1	Unused	
1	C2	(S)-AL 8810	10023
1	C3	15-deoxy-Δ ^{12,14} -Prostaglandin A ₂	10265
1	C4	3-methoxy Prostaglandin F _{1α}	15170
1	C5	Prostaglandin F _{2α} 1,9-lactone	16130
1	C6	11β-Prostaglandin F _{2α}	16520
1	C7	Carbaprostacyclin	18210
1	C8	Thromboxane B ₂	19030
1	C9	(+)-15-epi Cloprostenol	10006692
1	C10	CAY10509	10009167
1	C11	Unused	
1	C12	Unused	

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 06/16/2015

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

Product Information

Plate	Well	Contents	Item Number
1	D1	<i>Unused</i>	
1	D2	15-deoxy- $\Delta^{12,14}$ -Prostaglandin A ₁	10065
1	D3	15(R)-15-methyl Prostaglandin A ₂	10270
1	D4	11-deoxy Prostaglandin F _{1β}	15510
1	D5	5- <i>trans</i> Prostaglandin F _{2β}	16250
1	D6	AL 8810*	16735
1	D7	5- <i>cis</i> Carbaprostacyclin	18211
1	D8	Fluprostenol Lactone Diol	70037
1	D9	<i>ent</i> -Prostaglandin F _{2α}	10008122
1	D10	CAY10510	10009168
1	D11	<i>Unused</i>	
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	16,16-dimethyl Prostaglandin A ₁	10080
1	E3	16,16-dimethyl Prostaglandin A ₂	10280
1	E4	11-deoxy Prostaglandin F _{1α}	15512
1	E5	U-44069	16440
1	E6	Cloprostenol (sodium salt)	16764
1	E7	Ciprostene (calcium salt)	18216
1	E8	Fluprostenol serinol amide	10004236
1	E9	AL 8810 methyl ester	10008370
1	E10	Fluprostenol methyl ester	10010151
1	E11	<i>Unused</i>	
1	E12	<i>Unused</i>	
1	F1	<i>Unused</i>	
1	F2	AL 8810 isopropyl ester	10113
1	F3	16-phenoxy tetranor Prostaglandin A ₂	10285
1	F4	13,14-dihydro Prostaglandin F _{1α}	15610
1	F5	5- <i>trans</i> U-44069	16442
1	F6	(+)-Cloprostenol (sodium salt)	16766
1	F7	Prostaglandin I ₂ (sodium salt)	18220
1	F8	17-phenyl trinor Prostaglandin F _{2α} serinol amide	10004237
1	F9	8- <i>iso</i> -17-phenyl trinor Prostaglandin F _{2α}	10008435
1	F10	15-keto-17-phenyl trinor Prostaglandin F _{2α} ethyl amide	10010405
1	F11	<i>Unused</i>	
1	F12	<i>Unused</i>	

*Sold under license from Alcon

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

Product Information

Page 3 of 3



Plate	Well	Contents	Item Number
1	G1	<i>Unused</i>	
1	G2	5- <i>trans</i> Latanoprost (free acid)	10129
1	G3	17-phenyl trinor Prostaglandin A ₂	10288
1	G4	15-keto Prostaglandin F _{1α}	15710
1	G5	U-46619	16450
1	G6	15-keto Fluprostenol	16785
1	G7	Piriprost (potassium salt)	18228
1	G8	17-phenyl trinor Prostaglandin E ₂ serinol amide	10004238
1	G9	8- <i>iso</i> -17-phenyl trinor Prostaglandin F _{2β}	10008436
1	G10	AL 8810 ethyl amide	10010470
1	G11	<i>Unused</i>	
1	G12	<i>Unused</i>	
1	H1	<i>Unused</i>	
1	H2	Treprostinil	10162
1	H3	17-phenyl trinor-13,14-dihydro Prostaglandin A ₂	10290
1	H4	17,20-dimethyl Prostaglandin F _{1α}	15790
1	H5	5- <i>trans</i> U-46619	16452
1	H6	CAY10580	16835
1	H7	Beraprost (sodium salt)	18230
1	H8	(+)-5- <i>trans</i> Cloprostenol	10004970
1	H9	15(R)-Pinane Thromboxane A ₂	10008510
1	H10	17-phenoxy trinor Prostaglandin F _{2α}	10010839
1	H11	<i>Unused</i>	
1	H12	<i>Unused</i>	

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com